

## Product Information

### BTN-SRPK1

 Product Number : **04-460-20N**

#### Product description

Full-length human SRPK1 [1-655(end) amino acids and V504 deletion of accession number NP\_003128.3] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (78 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

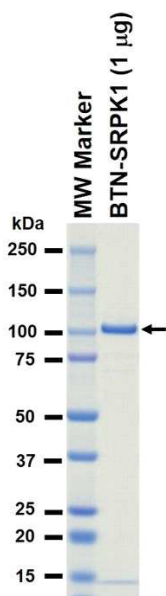
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
 1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.  
 Avoid repeating freeze-thaws.

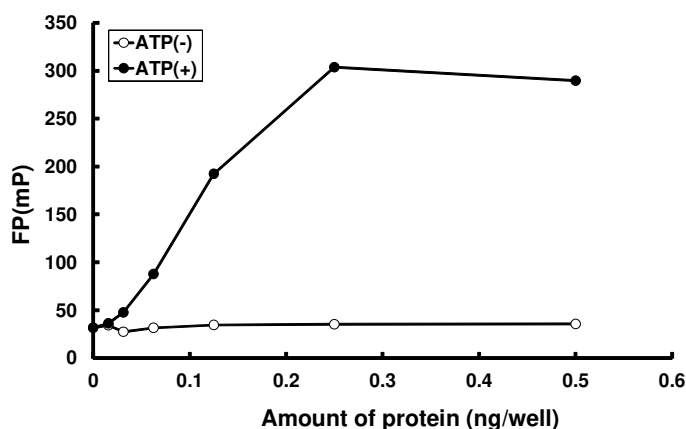
#### SDS-PAGE



Purity: 84 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : SRPKtide

 ATP : 100  $\mu$ M